

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit
Lot Number: K0082495
Expiration Date: 30 JUN 2026
Catalog Number: K156QO-Series, K151QO-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- γ Antibody	D0082162	2-8°C	30 APR 2027
Human IFN- γ Plate	Z0047968	2-8°C	30 SEP 2026
Proinflammatory Panel 1 (human) Calibrator Blend	A0080397	2-8°C	31 MAY 2029
Diluent 2	M0020129	$\leq -10^\circ\text{C}$	30 JUN 2026
Diluent 3	M0030127	$\leq -10^\circ\text{C}$	30 SEP 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	1400	pg/mL

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 JUN 2024